

XFD488 Anti-human CD22 Antibody *HIB22*

Catalog Number: 10220150, 10220151

Unit Size: 100 tests, 500 tests

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD22 (BL-CAM, Siglec-2)
Clone	HIB22
Conjugate	AF488

Biological Properties

Appearance	Orange liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	AF488
Excitation Wavelength	499 nm
Emission Wavelength	520 nm

Applications

HIB22 is an anti-human monoclonal antibody that recognizes the CD22 antigen. CD22 (also known as BL-CAM or Lyb8) is a 120 - 130 kD glycoprotein that is located on the surface of cells like B cells and stem cells. CD22 plays a role in essential cellular pathways, namely, the negative regulation of B cell receptor signaling pathway. Moreover, in many organisms, it is a negative regulator of calcium-mediated signaling, suppresses immunoglobulin secretion and is involved in the negative regulation of B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands like PI3-kinase, PLCg1, p53/56lyn and CD45. CD22 is a fairly uncommon antibody target, with a little more than 4000 publications in the last decade. Even still, CD22 is vital to cell biology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to XFD488 (ex/em = 499/520 nm). XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 488 nm laser and 533/30 nm bandpass filter (for example, as in the BD Accuri™ C6 Plus).